

Monolith Protocol MO-P-068

TGT Dimerization

Shigella bacteria constitute the causative agent of bacillary dysentery, an acute inflammatory disease causing the death of more than one million humans per year. The enzyme tRNA-guanine transglycosylase (TGT) catalyzes the base exchange of the genetically encoded guanine at the wobble position of particular tRNAs by the modified base preQ1. Further modification of preQ1 and the final incorporation of queuine into the tRNAs lead to the development of the virulence phenotype of *Shigella*, thus making TGT a suitable drug target. The enzyme is biologically active only as a homodimer, as one monomer subunit performs the catalytic base-exchange reaction while the second subunit is required to position the tRNA substrate in the correct orientation.

protein dimerization | homodimer

A1. Target/Fluorescent Molecule

Queuine tRNA-ribosyltransferase (TGT)

uniprot.org/uniprot/P28720

A2. Molecule Class/Organism

tRNA-ribosyltransferase

Zymomonas mobilis subsp. *mobilis* (strain ATCC 31821 / ZM4 / CP4)

A3. Sequence/Formula

GSMEATAQE TDRPRFSFSI AAREGKARTG TIEMKRGVIR TPAFMPVGTA ATVKALKPET VRATGADIIL GNTYHMLLRP
GAERIAKLGG LHSFMGWDRP ILTDSGGYQV MSLSSLTKQS EEGVTFKSHL DGSRHMLSPE RSIEIQHLLG SDIVMAFDEC
TPYPATPSRA ASSMERSMRW AKRSRDAFDS RKEQAENAAL FGIQQGSVFE NLRQQSADAL AEIGFDGYAV GGLAVGEGQD
EMFRVLDFSV PMLPDDKPHY LMGVGKPDDI VGAVERGIDM FDCVLPTRSG RNGQAFTWDG PINIRNARFS EDLKPLDSEC
HCAVCQKWSR AYIHXLRAG EILGAMLMT EILGAMLMT HNIIFYQQLM QKIRDSISEG RFSQFAQDFR ARYFARN S

wt X = H

H333D X = D

A4. Purification Strategy/Source

University of Marburg (Prof. Gerhard Klebe)¹

A5. Stock Concentration/Stock Buffer

wt

8.54 mg/ml | 199 µM

10 mM HEPES, pH 7.4, 1 M NaCl

H333D

7.76 mg/ml | 180 µM

10 mM HEPES, pH 7.4, 1 M NaCl

¹ Jakobi et al., ACS Chem. Biol., 10, 8, 1897-1907 (2015)

A6. Molecular Weight/Extinction Coefficient

<i>wt</i>	<i>H333D</i>
43,013.8 Da	42,991.8 Da
33,920 M ⁻¹ cm ⁻¹ (ε ₂₈₀)	33,920 M ⁻¹ cm ⁻¹ (ε ₂₈₀)

A7. Dilution Buffer

10 mM HEPES, pH 7.4, 1 M NaCl, 0.005% TWEEN® 20

A8. Labeling Strategy

Monolith Protein Labeling Kit RED – NHS 2nd Generation (MO-L011, NanoTemper Technologies GmbH)
1* Dye RED-NHS 2nd Generation (10 µg) | 1* B-Column

A9. Labeling Procedure

1. Prepare 100 µL of a 10 µL TGT solution in dilution buffer.
2. Add 25 µL of DMSO to 10 µg RED-NHS 2nd Generation dye to obtain a ~600 µM solution. Mix the dye thoroughly by vortexing and make sure that all dye is dissolved.
3. Mix 5 µL of the 600 µM dye solution with 95 µL of dilution buffer to obtain 100 µL of a 30 µM dye solution (3x protein concentration).
4. Mix TGT and dye in a 1:1 volume ratio (200 µL final volume, 2.5% final DMSO concentration).
5. Incubate for 30 minutes at room temperature in the dark.
6. In the meantime, remove the top cap of the B-Column and pour off the storage solution. Remove the bottom cap and place with adapter in a 15 mL tube.
7. Fill the column with dilution buffer and allow it to enter the packed resin bed completely by gravity flow. Discard the flow through collected. Repeat this step 3 more times.
8. Add 200 µL of the labeling reaction from step 4 to the center of the column and let sample enter the resin bed completely.
9. Add 500 µL of dilution buffer after the sample has entered and discard the flow through.
10. Place column in a new collection tube, add 500 µL of dilution buffer and collect the eluate.
11. Keep the labeled TGT (~2 µM) on ice in the dark.

A10. Labeling Efficiency

Measurement of protein concentration and degree of labeling (DOL) using a NanoDrop™:

nanotempertech.com/dol-calculator

<i>wt</i>			
Absorbance A ₂₀₅	1.66	Protein concentration	1.2 µM
Absorbance A ₆₅₀	0.05	Degree-of-labeling (DOL)	0.29
<i>H333D</i>			
Absorbance A ₂₀₅	1.54	Protein concentration	1.1 µM
Absorbance A ₆₅₀	0.07	Degree-of-labeling (DOL)	0.49

B1. Ligand/Non-Fluorescent Binding Partner

Queuine tRNA-ribosyltransferase (TGT)

uniprot.org/uniprot/P28720

B2. Molecule Class/Organism

tRNA-ribosyltransferase

Zymomonas mobilis subsp. *mobilis* (strain ATCC 31821 / ZM4 / CP4)

B3. Sequence/Formula

GSMVEATAQE TDRPRFSFSI AAREGKARTG TIEMKRGVIR TPAFMPVGTA ATVKALKPET VRATGADIIL GNTYHMLRP
GAERIAKLGG LHSFMGWDRP ILTDSGGYQV MSLSSLTKQS EEGVTFKSHL DGSRHMLSPE RSIEIQHLLG SDIVMAFDEC
TPYPATPSRA ASSMERSMRW AKRSRDAFDS RKEQAENAAL FGIQQGSVFE NLRQQSADAL AEIGFDGYAV GGLAVGEGQD
EMFRVLDFSV PMLPDDKPHY LMGVGKPDII VGAVERGIDM FDCVLPTRSG RNGQAFTWDG PINIRNARFS EDLKPLDSEC
HCAVCQKWSR AYIHXLIRAG EILGAMLMT EHNIAFYQQLM QKIRDSISEG RFSQFAQDFR ARYFARN

wt X = H

H333D X = D

B4. Purification Strategy/Source

University of Marburg (Prof. Gerhard Klebe)²

B5. Stock Concentration/Stock Buffer

wt

8.54 mg/ml | 199 μ M

10 mM HEPES, pH 7.4, 1 M NaCl

H333D

7.76 mg/ml | 180 μ M

10 mM HEPES, pH 7.4, 1 M NaCl

B6. Molecular Weight/Extinction Coefficient

wt

43,013.8 Da

33,920 M⁻¹cm⁻¹ (ϵ_{280})

H333D

42,991.8 Da

33,920 M⁻¹cm⁻¹ (ϵ_{280})

B7. Serial Dilution Preparation

1. Prepare a PCR-rack with 16 PCR tubes. Add 20 μ L of 40 μ M TGT into tube **1**. Then, transfer 10 μ L of dilution buffer into tubes **2** to **16**.
2. Prepare a 1:1 serial dilution by transferring 10 μ L from tube to tube. Mix carefully by pipetting up and down. Remember to discard 10 μ L from tube **16** to get an equal volume of 10 μ L for all samples.
3. Mix 2 μ L of labeled TGT (~2 μ M) with 198 μ L of dilution buffer to obtain 200 μ L of ~20 nM TGT.
4. Add 10 μ L of labeled TGT (~20 nM) to each tube from **16** to **1** and mix by pipetting.
5. Incubate tubes for at least 30 minutes³ at room temperature in the dark before loading capillaries.

² Jakobi et al., *ACS Chem. Biol.*, 10, 8, 1897-1907 (2015)

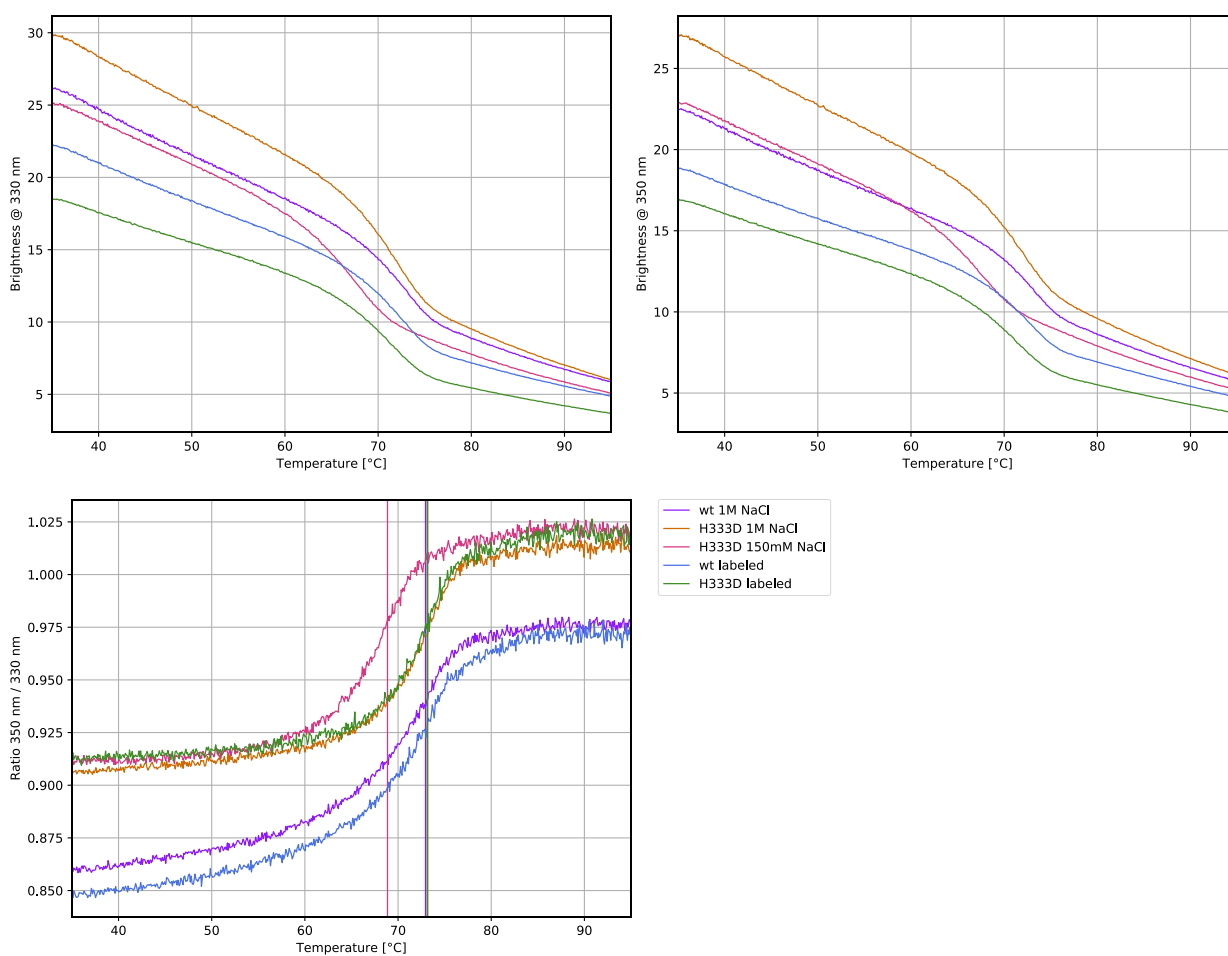
³ As the decay of stable dimers can be very slow, longer incubation times may be required to reach complete equilibrium.

C. Applied Quality Checks

Validation of structural integrity of unlabeled and labeled TGT in low and high salt buffers using Tycho NT.6:

nanotempertech.com/tycho

wt 1M NaCl	10 μ L of 2 μ M wt in dilution buffer	$T_i = 72.9^\circ\text{C}$
H333D 1M NaCl	10 μ L of 2 μ M H333D in dilution buffer	$T_i = 73.2^\circ\text{C}$
H333D 150 mM NaCl	10 μ L of 2 μ M H333D in 10 mM HEPES, pH 7.4, 150 mM NaCl	$T_i = 68.9^\circ\text{C}$
wt labeled	10 μ L of wt B-column eluate ($\sim 2 \mu\text{M}$)	$T_i = 73.1^\circ\text{C}$
H333D labeled	10 μ L of H333D B-column eluate ($\sim 2 \mu\text{M}$)	$T_i = 73.2^\circ\text{C}$



D1. MST System/Capillaries

Monolith NT.115^{Pico} Red (NanoTemper Technologies GmbH)

Premium Capillaries Monolith NT.115 (MO-K025, NanoTemper Technologies GmbH)

D2. MST Software

MO.Control v1.6 (NanoTemper Technologies GmbH)

nanotempertech.com/monolith-mo-control-software

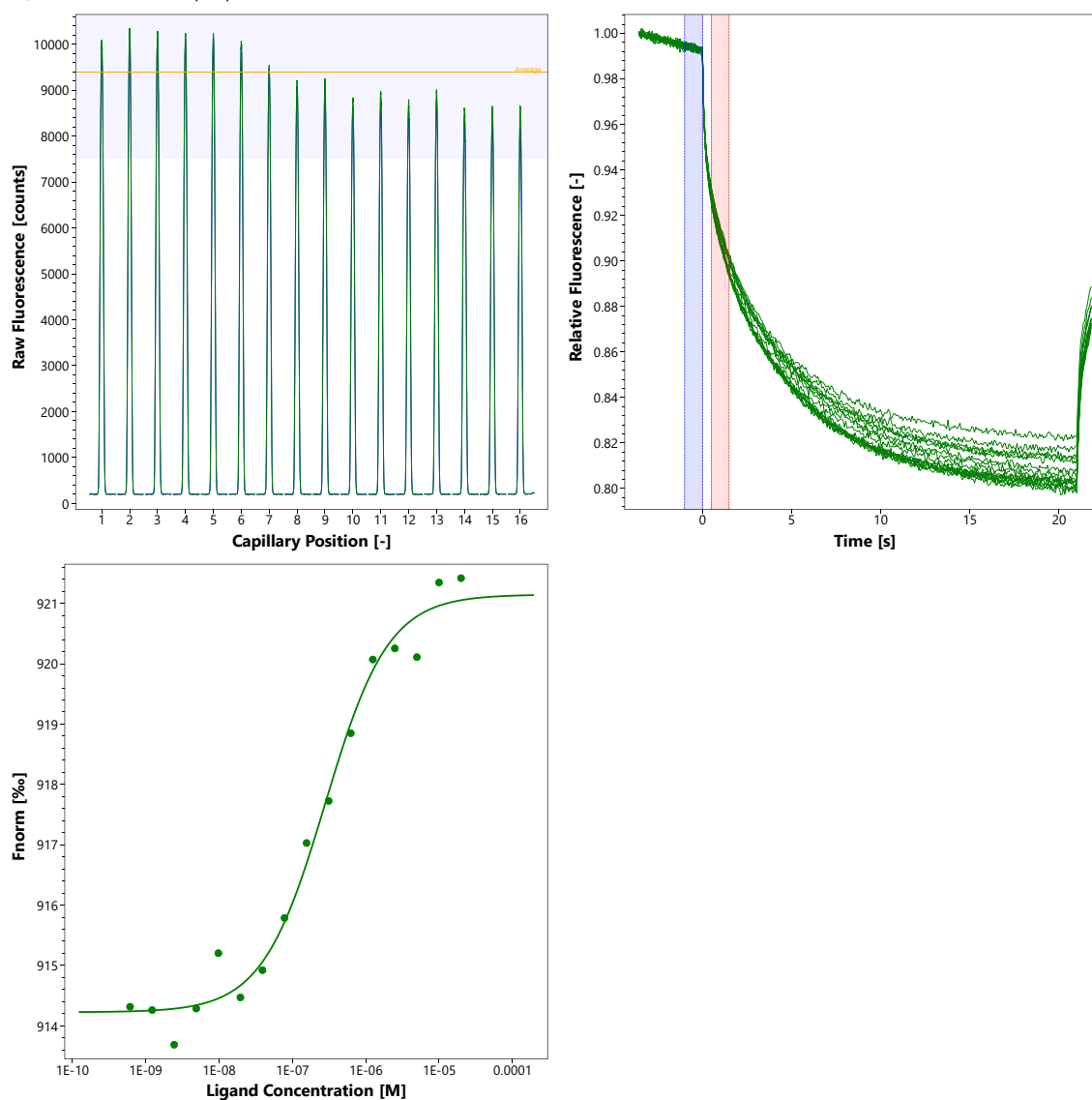
D3. MST Experiment (Assay Buffer/Concentrations/Temperature/MST Power/Excitation Power)

10 mM HEPES, pH 7.4, 1 M NaCl, 0.005% TWEEN® 20

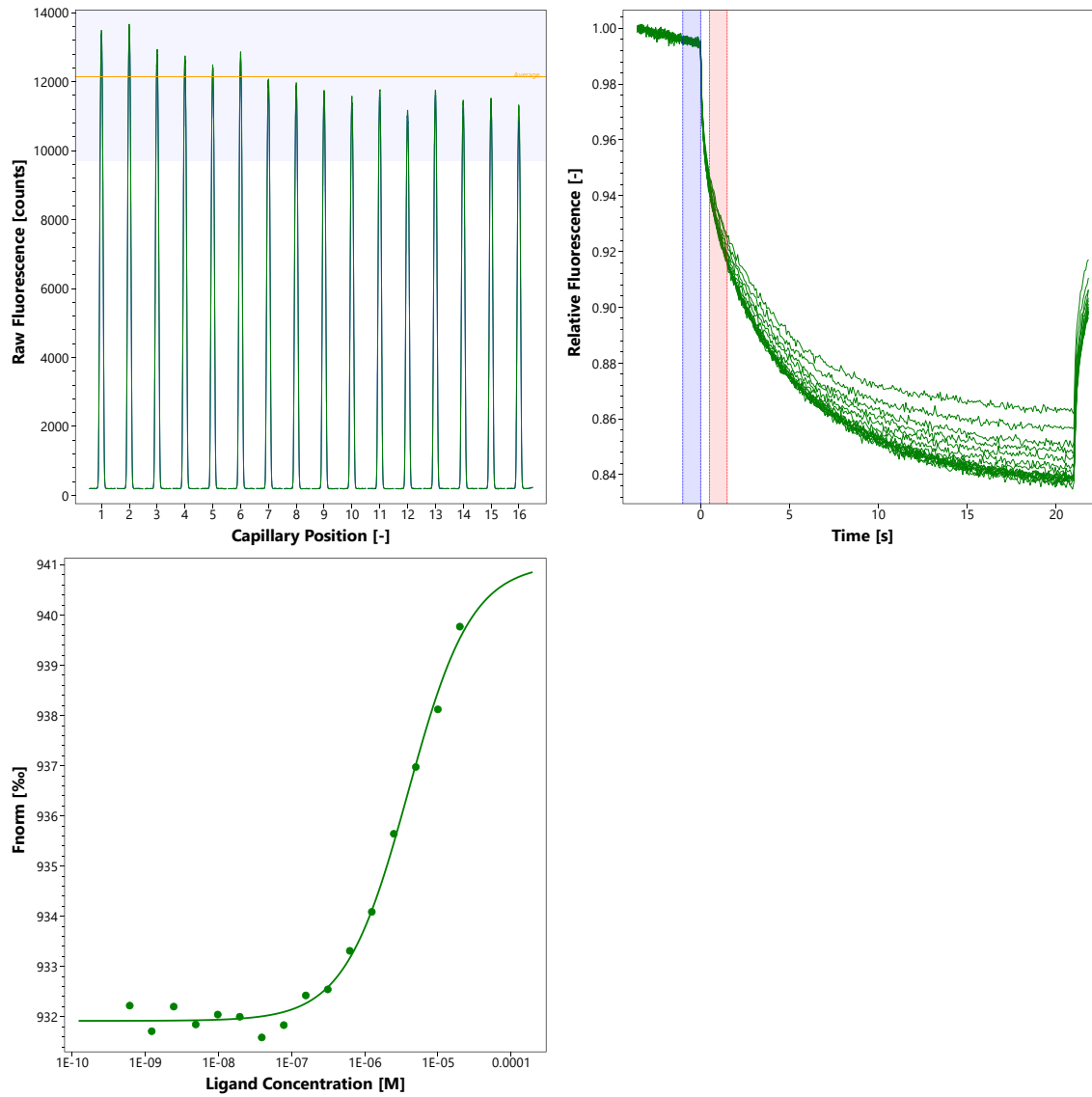
10 nM (labeled) TGT | 20 µM – 0.6 nM (unlabeled) TGT | 24°C | medium MST power | 20% excitation power

D4. MST Results (Capillary Scan/Time Traces/Dose Response)

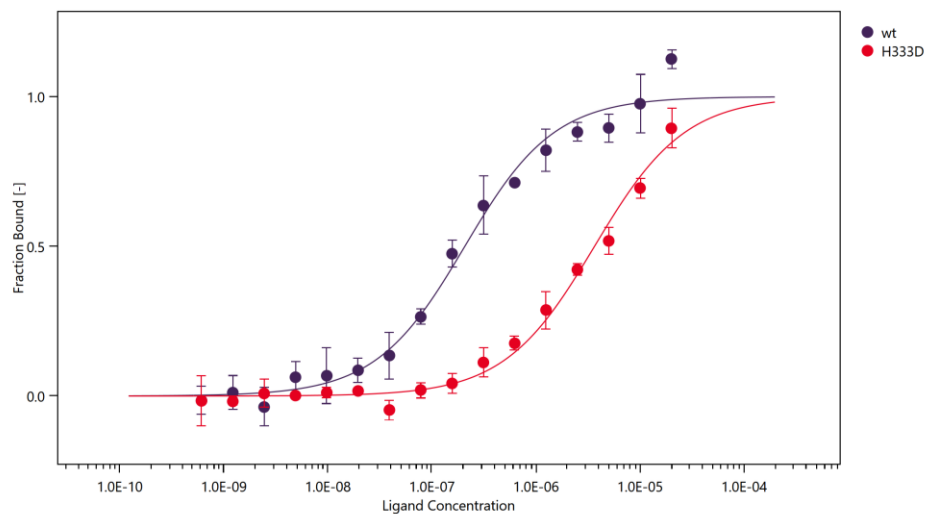
$K_d = 202 \pm 33$ nM (wt)



$K_d = 3.6 \pm 0.6 \mu\text{M}$ (H333D)



Comparison of dimerization affinities wt vs. H333D (n = 3):



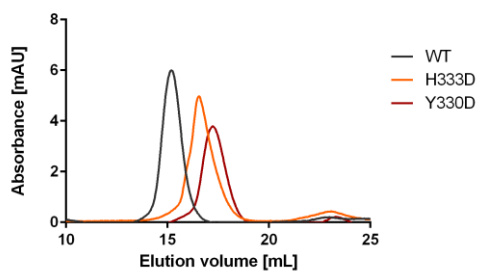
D5. Reference Results/Supporting Results

10 μ M wt: 2% monomer / 98% dimer
 10 μ M H333D: 58% monomer / 42% dimer
 10 μ M Y330D: 99% monomer / 1% dimer

Native MS

[Jakobi et al., ACS Chem. Biol. 2015, 10, 8, 1897-1907 \(2015\)](#)

Analytical SEC:



E. Contributors

Dzung Nguyen⁴, Andreas Langer⁵

⁴ Institute of Pharmaceutical Chemistry (IPC), University of Marburg, Marburg, Germany

⁵ NanoTemper Technologies GmbH, München, Germany | nanotempertech.com